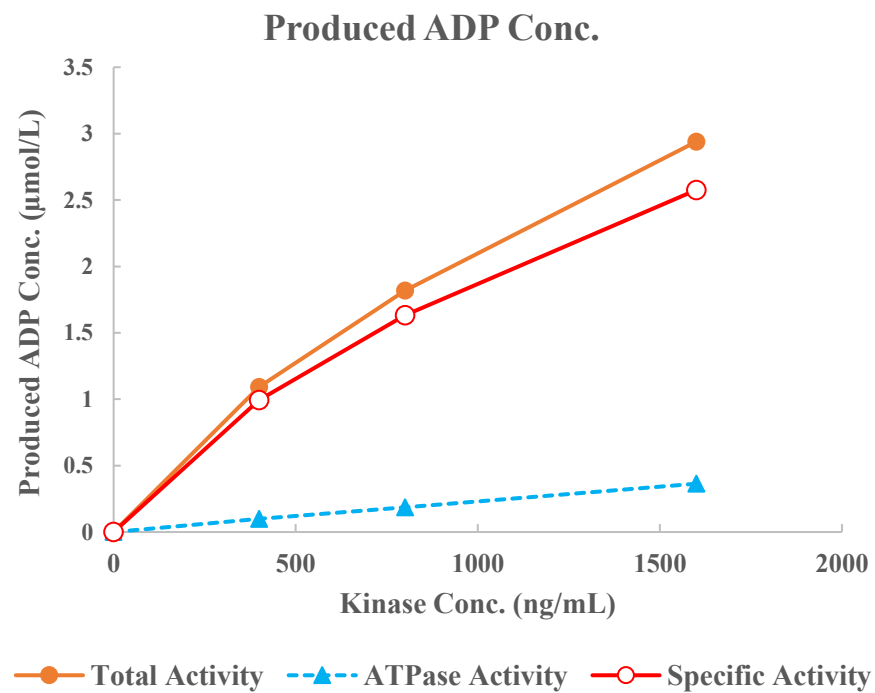
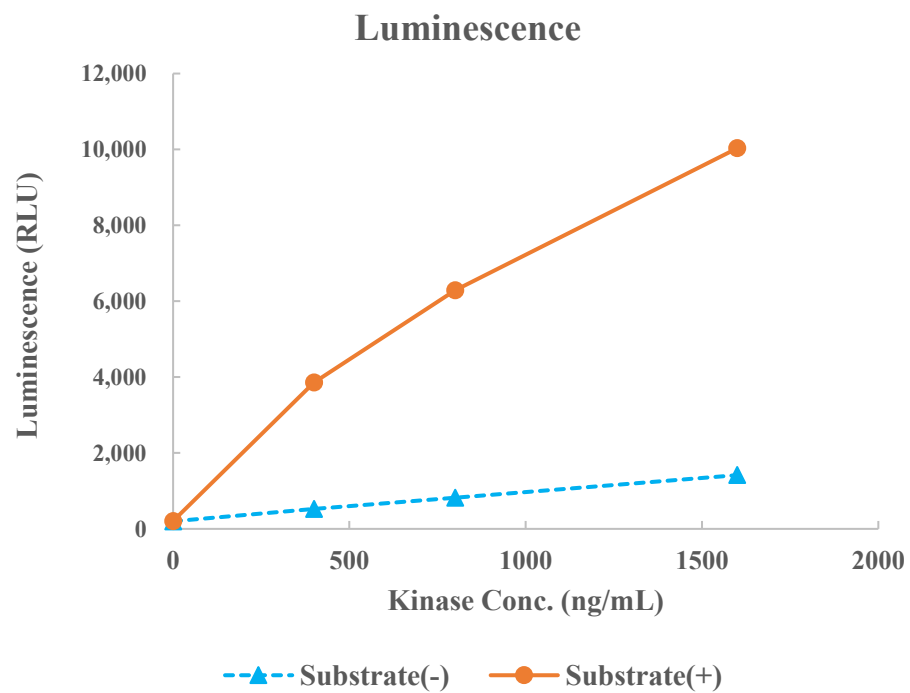


p70S6K β (RPS6KB2)



Kinase	p70S6K β (RPS6KB2), 01-155
Substrate	Biotin-S6K2 peptide, BTSU14
Substrate Conc.	10 μ mol/L
ATP Conc.	5 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour